

Version 2b Last updated 4 September 2020

ab270048

His-Tag Protein Expression Check Kit

A product of Expedeon, an
Abcam company

Applicable to Expedeon product codes 4003-0030.

View ab270048

His-Tag Protein Expression Check Kit:

www.abcam.com/ab270048

(use www.abcam.cn/ab270048 for China, or <http://www.abcam.co.jp/ab270048> for Japan)

For the rapid qualitative detection of His-tagged proteins.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	2
2. Materials Supplied and Storage	4
3. Technical Considerations	5
4. Assay Procedure	7

1. Overview

His-Tag Protein Expression Check Kit (ab270048) is a quick competitive immunochromatography test that allows users to verify and monitor successful expression of His-tagged recombinant proteins before undertaking purification, as well as identifying fractions of most interest before performing other more labor intensive techniques such as western blot or ELISA.

The assay provides a qualitative yes/no answer, although a semi-quantitative measurement of the protein abundance in the sample can be determined using internal references and controls.

A His-tagged protein on the “Test line” (T line) and a ULFA-tagged protein on the “Control line” (C line) are immobilized on the nitrocellulose strips.

40 nm Gold conjugated anti-HisTag antibody and 40 nm Gold conjugated anti-ULFA tag conjugates are dried on the conjugate pad and will selectively bind to the T and C line, respectively. The C line is assay-independent and should always appear as a strong red line; if it is not visible, the test is not valid and should be repeated.

His-Tag Protein Expression Check Kit (ab270048) is formulated as a competitive assay so that when the sample does not contain any His-tagged protein, the T line will appear as a strong red line (see Figure 1-A). On the contrary, the T line will not be visible (or appear as a pale red line) if the sample contains the His- tagged protein that competes with the binding of the Gold conjugated anti-HisTag antibody conjugate to the antigen immobilized on the T line (see Figure 1-B).

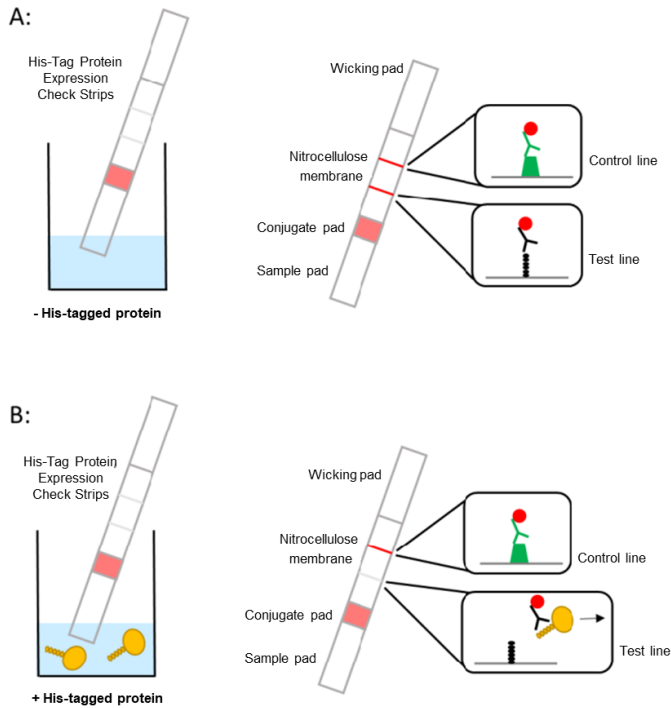


Figure 1. His-Tag Protein Expression Check Kit (ab270048) assay procedure. A: results in the absence of His-tagged protein, B: results in presence of His-tagged proteins.

2. Materials Supplied and Storage

Upon receipt, store kits at +4°C. Kit can be stored for 1 year from receipt, if components have not been reconstituted.

Avoid repeated freeze-thaws of reagents.

Item	30 tests	Storage temperature
10X Running Buffer	1 vials	+4°C
His-Tag Protein Expression Check Strips	30 Strips	+4°C

Not supplied: 96-wells low binding plate.

3. Technical Considerations

3.1 Sample Considerations:

The His-Tag Protein Expression Check Kit (ab270048) has been optimized with proteins containing accessible consecutive histidine regions at the amino- or carboxyl- terminus exposed in their native/ assay conformation.

The strips are compatible with cell culture media and lysate, as well as most common components used in purification processes (see table below).

Assay compatibility and interfering substances:

Substance	Compatibility
1X PBS	Fully compatible
1X TBS	Fully compatible
RIPA buffer	Fully compatible
Tween20	< 5%
CHAPS	< 1%
Triton	< 5%
SDS	< 1%
Imidazole	< 125 mM
Guanidine HCl	< 100 mM
Urea	< 125 mM
DTT	< 100 mM
2-mercaptoethanol	< 100 mM
EDTA	< 100 mM

It is advisable to run the samples in duplicate.

A sample without the His-tagged protein should always be added as negative control: the strip will show the highest signal intensity on the T line that can be expected. The C line should always have the same signal intensity. To confirm the presence of expressed His-

tagged protein in the sample, a decrease in signal intensity of the T line should be seen, as compared to the negative control sample. Reference samples of known His-tagged protein concentration can be used for a semi-quantitative analysis.

3.2 Assay Considerations:

- The strips are single-use.
- Always store the unused strips in the closed desiccant pot to prevent moisture from compromising their functionality.
- Make sure the flow is consistent and that both the sample pad overlapping the conjugate pad, and the conjugate pad overlapping the nitrocellulose membrane are making physical contact. If not, a slight bend of the strip (avoiding touching or damaging the nitrocellulose membrane) is enough to restore the contact and establish a steady and effective flow.

4. Assay Procedure

- 4.1 Bring all the kit components to room temperature.
- 4.2 Dilute the 10X Running Buffer with distilled water down to 1X.
- 4.3 Only 80 μ L of diluted sample are required for each strip. Dilute the lysate/sample containing the His-tagged protein down to 0.01-0.5 mg/mL in 1X Running Buffer; the appropriate range is protein dependent and should be determined empirically. If the protein concentration in the sample is unknown, perform serial 1:2 or 1:3 dilutions in 1X Running Buffer.
- 4.4 Load 80 μ L of diluted protein into a 96-well clear plate (low protein binding) and dip the end of the strip with the sample pad into the liquid.
- 4.5 Wait 10-15 minutes for the T and C lines to develop (do not let the strips to dry out before checking the result).
- 4.6 Check the strips; the intensity of the T lines inversely correlates with the amount of His tagged protein present in the sample.

Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)